



CERTIFICATE OF ANALYSIS

COA – 04417

Product Description: MSD® Performance Qualification Kit, PR

Catalog #/Size: R21QQ-2

Lot Number: PQ100020

Expiration Date: 30 APR 2015

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD’s specifications.

SECTOR PR2 Plate Lot Number: 01D01A

SECTOR PR2 Plate Catalog Number: L13XA

SECTOR PR2 Plate Expiration Date: 30 APR 2015

Free Tag 15 000 Lot Number: Y0260146

Free Tag 15 000 Part Number: N/A

Free Tag 15 000 Expiration Date: 31 JUL 2017

Read Buffer T(1x) Lot Number: Y0110053

Read Buffer T(1x) Catalog Number: R92TA

Read Buffer T(1x) Expiration Date: 31 JAN 2017

Specifications:

	Max	Min
Free Tag Mean	18 000	12 000
Free Tag CV %	6.0	0
Dark Signal Mean	16	-16
Dark Noise (SD)	40	0

MSD QC Results: Using instrument: PR400

S/N: 1150040915035

Free TAG Mean: 13 154

Dark Signal Mean: -1

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		04 SEP 2014
Review/Approval	Anastasia Belilovets	Quality Control		04 SEP 2014
Review/Approval	Richard Mann	Quality Assurance		04 SEP 2014